

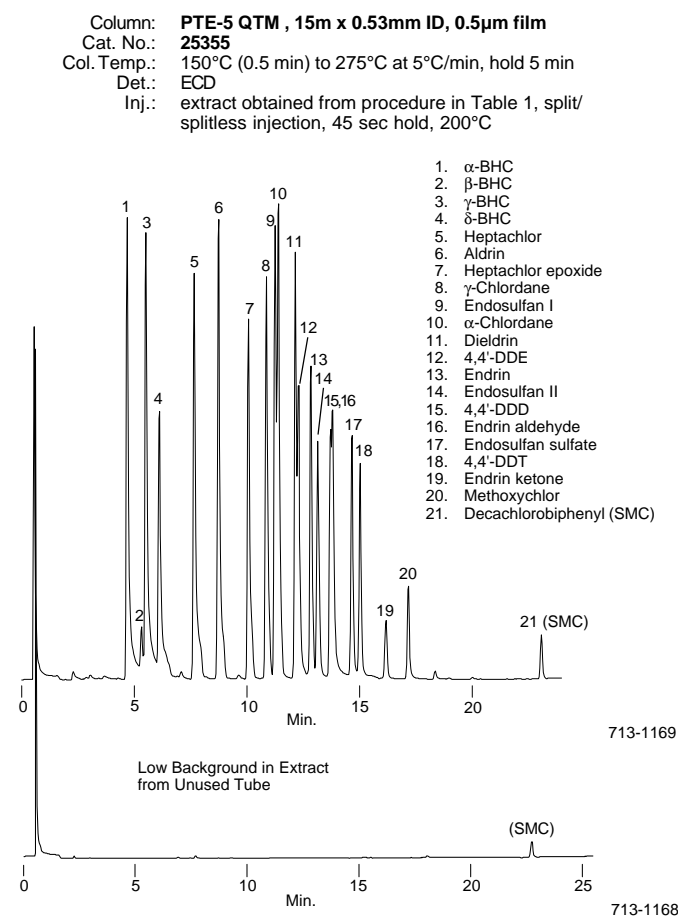
Low Background Solid Phase Extraction and Capillary GC Analysis of Chlorinated Pesticides

Solid phase extraction has been the main focus for sample preparation for monitoring semivolatiles, including pesticides, in aqueous hazardous waste samples because it provides rapid, on-site sample preparation and analysis. Supelco's ENVI-8 polymerically bonded octylsilane-silica packing provides good performance for solid phase extraction of pesticides from aqueous samples.

Key Words:

- chlorinated pesticides • solid phase extraction
- aqueous hazardous waste • ENVI-8 silica packing

Figure A. Extracted Pesticides
(Low Concentration, 10ng each/100mL sample)



Since the early 1980s, the US Environmental Protection Agency's Office of Solid Waste has been formulating new methods for identifying potentially hazardous abandoned waste sites. Currently, the EPA is developing a series of "quick turnaround" (QTM) methods, intended for use as screening methods, to allow decisions to be made quickly about the extent and nature of contamination at a site.

In EPA investigations, an octylsilane-modified, silica-based packing material has shown promise for solid phase extraction (SPE) of pesticides. We have developed a polymerically bonded octylsilane-silica packing, ENVI™-8, which performs well for extracting pesticides from aqueous samples. The polymeric bonding allows the material to resist pH extremes that might be encountered in sampling, and the high carbon loading should enhance the packing's capacity for nonpolar analytes. Extractions performed in our laboratories will be described here.

Any screening method for pesticides will require an extremely inert sample preparation system, because these analytes are of concern at very low concentrations – 10-250ng/100mL – and are monitored by GC with electron capture detection. To ensure reliable extractions, we used ENVI-8 packing in glass tubes with Teflon® frits. Extractions were performed using a Visiprep™ DL Solid Phase Extraction Vacuum Manifold[®] (fitted with disposable Teflon sample guides) and a Visidry™ Drying Attachment[®]. The samples were introduced directly into the tubes using custom glass tube adapters and the Visiprep large volume sampler.

Table 1. Solid Phase Extraction of Chlorinated Pesticides from Aqueous Hazardous Wastes

Extraction Tube:	0.5g ENVI-8 packing in 6mL glass tube, Teflon frits
Conditioning:	3mL methanol 2mL 5% methanol in water
Sample:	100mL aqueous sample, pH adjusted to 5.0-7.0
Addition:	add 5mL methanol add 10µL of 5.0µg/mL decachlorobiphenyl in methanol (SMC) introduce using Visiprep large volume sampler pass through SPE tube at 5mL/min.
Drying:	dry tubes under nitrogen flow for 2-3 minutes, using Visidry Drying Attachment
Elution:	add 5mL hexane:acetone (90:10), allow to soak into packing bed allow 2 minutes static extraction (no vacuum) add 5mL hexane:acetone (90:10), begin dropwise elution (low vacuum) reduce sample volume to 1mL, using nitrogen

Table 2. High Recovery of Pesticides by Solid Phase Extraction

Pesticide	SPE Packing Capacity Level (250ng each pesticide)			Contamination Level (10ng each pesticide)		
	% Recovery ^a	Std. Dev. ^b	Coef. Variat.	% Recovery ^a	Std. Dev. ^b	Coef. Variat.
α-BHC	91.4	7.0	7.7	84.6	9.2	10.9
β-BHC	100.1	10.1	10.1	— ^c	—	—
γ-BHC	94.0	6.1	6.4	88.2	10.8	12.2
δ-BHC	92.9	7.2	7.7	85.1	7.3	8.5
Heptachlor	95.0	5.7	6.0	86.6	6.4	7.3
Aldrin	94.9	4.8	5.1	84.7	13.9	16.4
Heptachlor epoxide	96.8	4.2	4.4	96.3	5.6	5.8
γ-Chlordane	97.4	4.5	4.6	85.8	6.3	7.3
Endosulfan I	94.6	4.7	4.9	84.8	8.5	10.1
α-Chlordane	96.4	4.6	4.7	90.2	7.4	8.2
Dieldrin	98.2	3.8	3.9	88.6	6.0	6.8
4,4'-DDE	99.9	4.2	4.2	80.1	7.7	9.6
Endrin	97.9	13.2	13.5	77.8	20.1	25.9
Endosulfan II	98.2	3.8	3.8	90.0	8.0	8.9
4,4'-DDD	104.4	5.4	5.2	98.6 ^d	10.9	11.1
Endrin aldehyde	101.0	5.5	5.5	98.6 ^d	10.9	11.1
Endosulfan sulfate	104.6	3.9	3.7	96.2	7.9	8.3
4,4'-DDT	107.9	4.9	4.5	95.3	13.4	14.1
Endrin ketone	110.4	14.3	12.9	104.6	29.9	28.6
Methoxychlor	108.4	3.2	2.9	93.4	13.1	14.0
Decachlorobiphenyl (SMC)	101.0	7.0	6.9	72.9	14.0	19.3

^a Mean for 5 lots.^b n = 10 extractions.^c Could not be monitored under the GC conditions used (split-splitless, heated injection); cold on-column injection should reduce problem.^d Compounds coelute, area counts not separable at 10ng/mL each.

We packed 6mL glass tubes with 0.5g of ENVI-8 packing (retained with Teflon frits) and prepared pesticide standards for extraction, spiking the samples at 250ng each pesticide/100mL ("packing capacity level") and 10ng each pesticide/100mL ("contamination level"). The higher level would test whether the SPE tubes have adequate capacity for the analytes and the lower level would confirm that background levels are lower than the minimum detection levels for the analytes. After monitoring the effects of parameters such as pH and drying time on the extraction, we found the extraction procedure described in Table 1 to be excellent for this purpose.

Using the procedure in Table 1, we obtained high recovery levels for pesticides at both the high concentration and low concentration levels (Table 2). In general, recovery rates were very satisfactory, although they were somewhat variable for endrin and endrin ketone. Overall, the SPE tubes and the extraction procedure provided good results which met the criteria needed for pesticides extractions. Low contaminant level extractions exhibited few interferences, and lot-to-lot variability in recovery rates was acceptable.

If you are interested in examining the potential of these highly inert ENVI-8 solid phase extraction tubes for pesticides or other sample preparation procedures, please refer to our catalog or call our technical service chemists. You can also get information about other SPE tubes for environmental samples, and for descriptions of our SPE processing apparatus.

Ordering Information:

Description	Cat. No.
ENVI-8 SPE Tubes	
6mL glass tube, 0.5g packing, pk. of 30 Teflon frits	57107
Visiprep DL Disposable Liner SPE Vacuum Manifold	
12-port model	57044
24-port model	57265
Disposable Teflon Solvent Guides	
package of 100	57059
Visidry Drying Attachment	57100-U
Visiprep Large Volume Sampler	
for 4 samples	57275
Adapters	
for glass tubes	custom
PTE™-5 QTM Capillary Column	
1.5m x 0.53mm ID, 0.5µm phase film	25355

TrademarksENVI, PTE, Visidry, Visiprep – Sigma-Aldrich Co.
Teflon – E.I. du Pont de Nemours & Co., Inc.

For pesticide standards, refer to the Supelco general catalog.

•US Pat. Nos. D.289, 861; 4,810,471; other patents pending.

•US Pat. No. 4,810,471; other patents pending.

Note 64

For more information, or current prices, contact your nearest Supelco subsidiary listed below. To obtain further contact information, visit our website (www.sigma-aldrich.com), see the Supelco catalog, or contact Supelco, Bellefonte, PA 16823-0048 USA.

ARGENTINA • Sigma-Aldrich de Argentina, S.A. • Buenos Aires 1119 **AUSTRALIA** • Sigma-Aldrich Pty. Ltd. • Castle Hill NSW 2154 **AUSTRIA** • Sigma-Aldrich Handels GmbH • A-1110 Wien
BELGIUM • Sigma-Aldrich N.V./S.A. • B-2880 Bornem **BRAZIL** • Sigma-Aldrich Quimica Brasil Ltda. • 01239-010 São Paulo, SP **CANADA** • Sigma-Aldrich Canada, Ltd. • 2149 Winston Park Dr., Oakville, ON L6H 6J8
CZECH REPUBLIC • Sigma-Aldrich s.r.o. • 186 00 Praha 8 **DENMARK** • Sigma-Aldrich Denmark A/S • DK-2665 Vallensbaek Strand **FINLAND** • Sigma-Aldrich Finland/YA-Kemia Oy • FIN-00700 Helsinki
FRANCE • Sigma-Aldrich Chimie • 38297 Saint-Quentin-Fallavier Cedex **GERMANY** • Sigma-Aldrich Chemie GmbH • D-82041 Deisenhofen **GREECE** • Sigma-Aldrich (o.m.) Ltd. • Ilioupoli 16346, Athens
HUNGARY • Sigma-Aldrich Kft. • H-1067 Budapest **INDIA** • Sigma-Aldrich Co. • Bangalore 560 048 **IRELAND** • Sigma-Aldrich Ireland Ltd. • Dublin 24 **ISRAEL** • Sigma Israel Chemicals Ltd. • Rehovot 76100
ITALY • Sigma-Aldrich s.r.l. • 20151 Milano **JAPAN** • Sigma-Aldrich Japan K.K. • Chuo-ku, Tokyo 103 **KOREA** • Sigma-Aldrich Korea • Seoul **MALAYSIA** • Sigma-Aldrich (M) Sdn. Bhd. • Selangor
MEXICO • Sigma-Aldrich Química S.A. de C.V. • 50200 Toluca **NETHERLANDS** • Sigma-Aldrich Chemie BV • 3330 AA Zwijndrecht **NORWAY** • Sigma-Aldrich Norway • Torshov • N-0401 Oslo
POLAND • Sigma-Aldrich Sp. z o.o. • 61-663 Poznań **PORTUGAL** • Sigma-Aldrich Química, S.A. • Sintra 2710 **RUSSIA** • Sigma-Aldrich Russia • Moscow 103062 **SINGAPORE** • Sigma-Aldrich Pte. Ltd.
SOUTH AFRICA • Sigma-Aldrich (pty) Ltd. • Jet Park 1459 **SPAIN** • Sigma-Aldrich Química, S.A. • 28100 Alcobendas, Madrid **SWEDEN** • Sigma-Aldrich Sweden AB • 135 70 Stockholm
SWITZERLAND • Supelco • CH-9471 Buchs **UNITED KINGDOM** • Sigma-Aldrich Company Ltd. • Poole, Dorset BH12 4QH
UNITED STATES • Supelco • Supelco Park • Bellefonte, PA 16823-0048 • Phone 800-247-6628 or 814-359-3441 • Fax 800-447-3044 or 814-359-3044 • email:supelco@sial.com

H

Supelco is a member of the Sigma-Aldrich family. Supelco products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product for a particular use. Additional terms and conditions may apply. Please see the reverse side of the invoice or packing slip.

AQZ